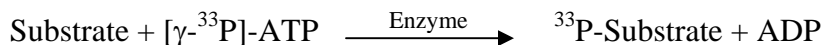


GRK6

(G protein-coupled receptor kinase 6, GPRK6)

- **CAT#:** GRK6
- **Enzyme:** Human GRK6
- **Substrate:** Casein, 1 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	46
Ro 31-8220	1,100
GW 5074	24,000
H-89	15,000

- **Scientific Information:** P43250 (<http://www.expasy.org/uniprot/P43250>)

Reference compound IC50 for GRK6

